

Product Number:	101736 (103159)
Lot Number:	501189 (80752)
Expiration:	16-Mar-2019
Storage Condition:	2-8°C

	PASS/FAIL
Release Criteria	Pass

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title


Signature/Date 29 MAR 2018

Statement: The above product was tested using Quanterix procedures and conforms to Quanterix's specifications.